

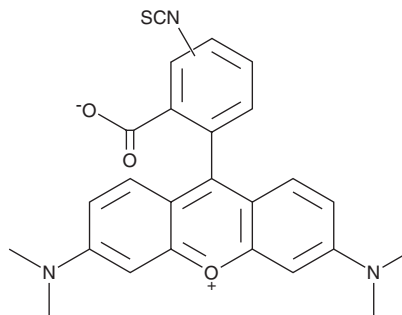
# PRODUCT INFORMATION



## Tetramethylrhodamine isothiocyanate (mixed isomers)

Item No. 19593

**CAS Registry No.:** 95197-95-8  
**Formal Name:** 9-(2-carboxyisothiocyanatophenyl)-3,6-bis(dimethylamino)-xanthylium, inner salt  
**Synonyms:** 5(6)-Tetramethylrhodamine isothiocyanate, 5(6)-TRITC, TRITC  
**MF:** C<sub>25</sub>H<sub>21</sub>N<sub>3</sub>O<sub>3</sub>S  
**FW:** 443.5  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 215, 256, 283, 356, 545 nm  
**Ex./Em. Max:** 552/575 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Tetramethylrhodamine isothiocyanate (TRITC) (mixed isomers) is supplied as a crystalline solid. A stock solution may be made by dissolving the TRITC (mixed isomers) in the solvent of choice, which should be purged with an inert gas. TRITC (mixed isomers) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of TRITC (mixed isomers) in these solvents is approximately 5 mg/ml.

TRITC (mixed isomers) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, TRITC (mixed isomers) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. TRITC (mixed isomers) has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

TRITC is an amine-reactive derivative of rhodamine that is used as a fluorescent label for antibodies and other probes (ex/em max = 552/575 nm).<sup>1</sup> It consists of a base tetramethylrhodamine molecule functionalized with an isothiocyanate reactive group at one of two hydrogen atoms on the bottom ring of the structure. TRITC is reactive towards primary amine groups on proteins, peptides, and other biomolecules.

### Reference

1. Li, B., Oestreich, S., and de Lange, T. Identification of human Rap1: Implications for telomere evolution. *Cell* **101**(5), 471-83 (2000).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/19/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
[WWW.CAYMANCHEM.COM](http://WWW.CAYMANCHEM.COM)